

ZytoLight®

FOXO1/PAX3 Dual Color Single Fusion Probe

For the detection of the translocation t(2;13)(q36;q14.1) by fluorescence *in situ* hybridization (FISH)

The ZytoLight® SPEC FOXO1/PAX3 Dual Color Single Fusion Probe is designed to detect the translocation t(2;13)(q36;q14.1) in alveolar rhabdomyosarcomas. Among solid tumors of the childhood, rhabdomyosarcoma is the most common soft tissue sarcoma.

Rhabdomyosarcomas are classified in two main categories: embryonal and alveolar rhabdomyosarcoma. The alveolar histology is associated with a poorer prognosis.

Ordering Information [Order Online »](#)

ZTV-Z-2018-50	50µl
ZTV-Z-2018-200	200µl

Manuals, SDS & CofA [View Online »](#)

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), using 10 µl probe solution per test.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).
For the Manual please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)